

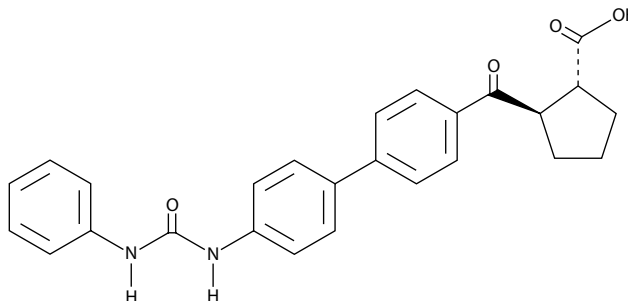
Product Information



A-922500

Item No. 10012708

CAS Registry No.: 959122-11-3
Formal Name: 2R-[[4'-[[[(phenylamino) carbonyl]amino][1R,1'-biphenyl]-4-yl-carbonyl]-cyclopentanecarboxylic acid
MF: C₂₆H₂₄N₂O₄
FW: 428.5
Purity: ≥95%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that A-922500 be stored as supplied at -20°C. It should be stable for at least one year. A-922500 is supplied as a crystalline solid. A stock solution may be made by dissolving the A-922500 in the solvent of choice. A-922500 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of A-922500 in ethanol is approximately 0.2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. A-922500 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, A-922500 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. A-922500 has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Diacylglycerol acyltransferases (DGATs) catalyze the final step in triglyceride synthesis. DGAT-1 deficient mice demonstrate a phenotype that is protective against the development of diet-induced obesity (DIO) or insulin resistance, implicating the actions of this enzyme in the development of such metabolic disorders. A-922500 is a potent orally active inhibitor of DGAT-1 activity, inhibiting both human and mouse forms of the enzymes with IC₅₀ values of 7 and 24 nM, respectively.¹ When administered at 3 mg/kg to DIO mice, A-922500 conferred significant weight loss within seven days without affecting food intake and significantly reduced plasma and liver triglycerides with chronic dosing.¹ A-922500 does not inhibit DGAT-2, ACAT-1, or ACAT-2.¹

Reference

1. Zhao, G., Souers, A.J., Voorbach, M., *et al.* Validation of diacyl glycerolacyltransferase I as a novel target for the treatment of obesity and dyslipidemia using a potent and selective small molecule inhibitor. *J. Med. Chem.* **51**, 380-383 (2008).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10012708

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/02/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com